

Product Notice

Description: VeriKine-HS™ Mouse Interferon Beta Serum ELISA Kit

Catalog No: 42410-1, 42410-2

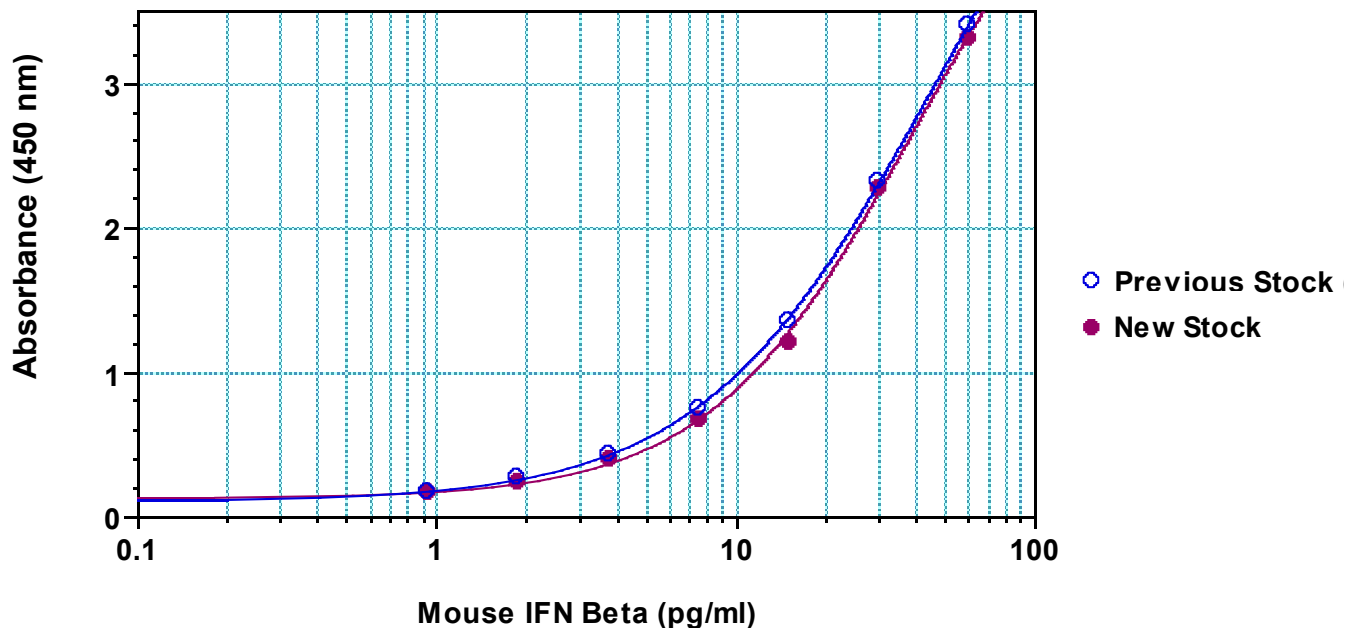
Lots: 7056 onward

Date: July 23, 2019

This notice is to inform you that a change has been made to the interferon standard used in the Mouse Interferon Beta ELISA Serum kit listed above. The formulation of the interferon standard remains the same, but the mouse interferon beta stock used in preparation of this standard has been replaced with a new stock. No changes to the remaining kit components have been implemented.

Please find example data below from a standard study comparing the new Mouse Interferon Beta stock with the previous stock in the 42410 ELISA kit.

IFN β Standard Stock Comparison in 42410



We thank you for your patience as PBL continues to improve the quality of its products. Please feel free to contact us at techservice@pbl assaysci.com if you have any product related questions.